

ProQinase™ BLK

BLK proto-oncogene, Src family tyrosine kinase

Recombinant Human Active Protein Kinase

HGNC Symbol: BLK

Synonyms: MODY11, MGC10442

Product No.: 0448-0000-1

Lot: 001

Description: Human BLK, full length, amino acids M1-P505 (as in [NCBI/Protein](#) entry NP_001706.2), N-terminal GST-HIS6 fusion protein with a Thrombin cleavage site, expressed in Sf9 insect cells

Product identity: BLK Lot 001, was confirmed as BLK by mass spectroscopy LC-ESI-MS/MS

Theoretical MW_{Fusion Protein}: 87,102 Da

Expression host: Sf9 insect cells

Purification: GST-Affinity Chromatography

Activation: This kinase was not activated by special procedures

Storage buffer: 50 mM TRIS-HCl pH 8.0, 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20 % glycerol

Storage temperature: -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10µl are not advisable. Avoid repeated freeze-thaw cycles!

Protein concentration: 0.120 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

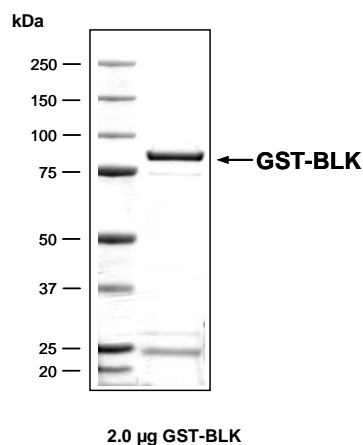
Biochemical Parameters:

Specific kinase activity (P_i transfer): 167 pmol/µg × min
ATP-K_M: 3.1 µM

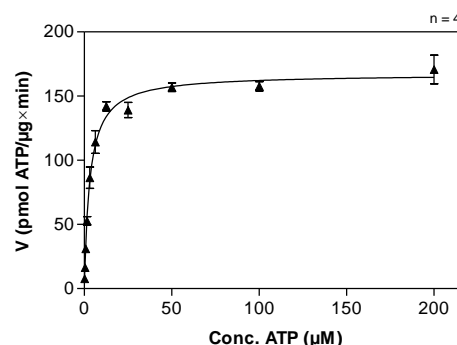
Additional assay technology:

BLK Lot 001 was also successfully tested by Reaction Biology for the use with the ADP-Glo™ Kinase assay from Promega ADP-Glo assay conditions may vary from radiometric assay conditions, please inquire for assay details

BLK Lot 001: Coomassie stain



BLK Lot 001: Determination of V_{max} and K_M value for ATP



- Assay conditions:
60 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 µM Na-orthovanadate
1.2 mM DTT
50 µg/ml PEG_{20,000}
ATP (variable)
Substrate: Poly(Glu:Tyr)_{4:1} 20 µg/ml
Kinase: 4 µg/ml
- Filter binding assay
MSFC membrane (Millipore)

ProQinase™ BLK

Product No.: 0448-0000-1

GST-BLK Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQGWQATF	GGGDHPPKSD	PMGHHHHHG	RRRASVAAGI	240
241	LVPRGSPGLD	GICSRMGLVS	SKKPDKEKPI	KEKDKGQWSP	LKVS AQDKDA	PPLPPLVFN	300
301	HLTPPPDEH	LDEDKHFVVA	LYDY TAMNDR	DLQMLKGEKL	QVLKGTGDWW	LARSLVTGRE	360
361	GYVPSNFVAR	VESLEMERWF	FRSQGRKEAE	RQLLAPINKA	GSFLIRESET	NKGAFSLSVK	420
421	DVTTQ GELIK	HYKIRCLDEG	GYIISPRITF	PSLQALVQHY	SKKGDGLCQR	LTLPCVRPAP	480
481	QNPWAQDEWE	IPRQSLRLVR	KLGSQGFGEV	WMGYKNNMK	VAIKTLKEGT	MSPEAFLGEA	540
541	NVMKALQHER	LVRLYAVVTK	EPIYIVTEYM	ARGCLLDFLK	TDEGSRLSLP	RLIDMSAQIA	600
601	EGMAYIERMN	SIHRDLRAAN	ILVSEALCCK	IADFGLARII	DSEYTAQEGA	KFPIKWTAPE	660
661	AIHFGVFTIK	ADVWSFGVLL	MEVVTYGRVP	YPGMSNPEVI	RNLERGYRMP	RPDTCPPELY	720
721	RGVIAECWRS	RPEERPTFEF	LQSVLEDFYT	ATERQYELQP			780

1-218: GST Red: HIS6-tag Pink: Thrombin cleavage site blue: BLK

BLK wt ¹ Amino Acid Sequence							
1	MGLVSSKKPD	KEKPIKEKDK	GQWSPLK VSA	QDKDAPPLPP	LVVFNHLTPP	PPDEHLDEDK	60
61	HFVVALYDYT	AMNDRDLQML	KGEKLQVLKG	TGDWWLARS L	VTGREGYVPS	NFVARVESLE	120
121	MERWFFRSQG	RKEAERQLLA	PINKAGSFLI	RESETNKGAF	SLSVKDVT TQ	GELIKHYKIR	180
181	CLDEGGYYIS	PRITFPSLQA	LVQHYSKKG D	GLCQRLTLPC	VRPAPQNPWA	QDEWEIPRQS	240
241	LRLVRKLGSG	QFGEVW MGY	KNNMKVAIKT	LKEGTMSPEA	FLGEANVMKA	LQHERLVRLY	300
301	AVVTKEPIYI	VTEYMARGCL	LDFLKTDEGS	RLSLPRLIDM	SAQIAEGMAY	IERMNSIHRD	360
361	LRAANILVSE	ALCCKIADFG	LARIIDSEYT	AQEGAKFPIK	WTAPEAIHFG	VFTIKADVWS	420
421	FGVLLMEVVT	YGRVPYPGMS	NPEVIRNLER	GYRMPRPDTC	PPELYRGVIA	ECWRSRPEER	480
481	PTFEFLQSVL	EDFYTATERQ	YELQP				540

blue: BLK sequence expressed in recombinant protein

¹[NCBI/Protein](#) accession number NP_001706.2